

Data Sheet

PDE7A

Murine, recombinant, N-terminal GST-tag

Catalog #: 60072

Lot#: 100916

Conc.: 0.16 mg/ml

Formulated in: 45 mM Tris-HCl, pH 8.0,
 124 mM NaCl, 2.4 mM KCl, 18 mM
 glutathione, 10% glycerol, and 3 mM DTT.

Stability: >6 months at -80°C

References:

1. Han, P., *et al.* (1998). *Genomics* Mar 1;**48(2)**:275-6.
2. Yang, G., *et al.* (2003). *J. Immunol.* Dec 15;**171(12)**:6414-20.
3. Yamamoto, *et al.* (2006) *Eur. J. Pharmacol.* Jul 10;**541(1-2)**:106-14.

Description:

Mouse PDE7A (GenBank Accession No. NM_001122759), a.a. 147-482(end), with N-terminal GST-tag, MW=65 kDa, expressed in a Baculovirus-infected Sf9 cell expression system.

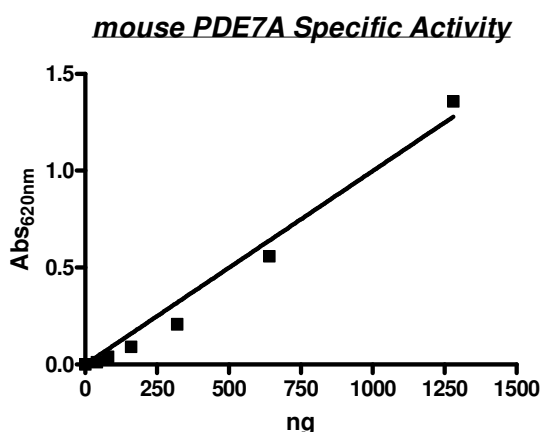
Specific Activity: 120 pmol/min/μg.

Assay condition: 10 mM Tris-HCl, pH 7.4, 10 mM MgCl₂, 0.05% Tween-20, 200 μM cAMP, 0.1 mg/ml BSA, and 0.8-26 ng/μl mouse PDE7A at 37°C for 20 min. Quantified by 5'-nucleotidase cleaving the 5'-AMP product and releasing the phosphate group which is detected by Malachite Green Reagent.

Application:

Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance



10% SDS-PAGE Coomassie staining

Lane 1:

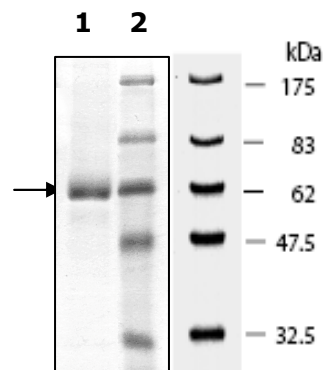
3 μg PDE7A

Lane 2:

Protein Marker
 BioLabs (#P7708L)

MW: 65 kDa.

Purity: >22%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com